

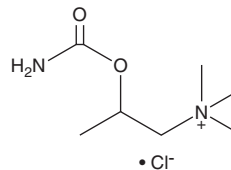
PRODUCT INFORMATION



Bethanechol (chloride)

Item No. 23830

CAS Registry No.: 590-63-6
Formal Name: 2-[(aminocarbonyl)oxy]-N,N,N-trimethyl-1-propanaminium, monochloride
Synonyms: (±)-Bethanechol, Carbamyl-β-methylcholine, NSC 30783
MF: C₇H₁₇N₂O₂ • Cl
FW: 196.7
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Bethanechol (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the bethanechol (chloride) in the solvent of choice. Bethanechol (chloride) is soluble in the organic solvent DMSO, which should be purged with an inert gas at a concentration of approximately 0.1 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of bethanechol (chloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of bethanechol (chloride) in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Bethanechol is an agonist of muscarinic acetylcholine receptors with IC₅₀ values of 1,837, 25, 631, 317, and 393 μM for M₁₋₅, respectively, in a radioligand binding assay using CHO cells expressing the human receptors.¹ It inhibits M₂-mediated increases in cyclic AMP induced by isoproterenol (Item No. 15592) in isolated guinea pig small intestine (IC₅₀ = 127 μM).² Bethanechol increases basal tone of isolated porcine intravesical ureter (EC₅₀ = 4.27 μM).³ It also induces fluid secretion in the ileum, duodenum, and jejunum of anesthetized rats when administered at a dose of 60 μg/kg.⁴ Formulations containing bethanechol have been used to increase urination and improve smooth muscle tone in the gastrointestinal tract.

References

1. Richards, M.H. and van Giersbergen, P.L. Human muscarinic receptors expressed in A9L and CHO cells: Activation by full and partial agonists. *Br. J. Pharmacol.* **114(6)**, 1241-1249 (1995).
2. Okamoto, H., Prestwich, S.A., Asai, S., *et al.* Muscarinic agonist potencies at three different effector systems linked to the M₂ or M₃ receptor in longitudinal smooth muscle of guinea-pig small intestine. *Br. J. Pharmacol.* **135(7)**, 1765-1775 (2002).
3. Hernández, M., Símonsén, U., Prieto, D., *et al.* Different muscarinic receptor subtypes mediating the phasic activity and basal tone of pig isolated intravesical ureter. *Br. J. Pharmacol.* **110(4)**, 1413-1420 (1993).
4. Young, A. and Levin, R.J. Intestinal hypersecretion of the refeed starved rat: A model for alimentary diarrhoea. *Gut* **33(8)**, 1050-1056 (1992).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/03/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM